

# SARTORIUS

## Simplifying Progress

## Diversity Fact Sheet

Sartorius 2022



---

“At Sartorius, diversity is not about statistics. It’s about culture. We foster diversity because it ultimately teaches us to be more open. And we consider openness as a key prerequisite for success.”

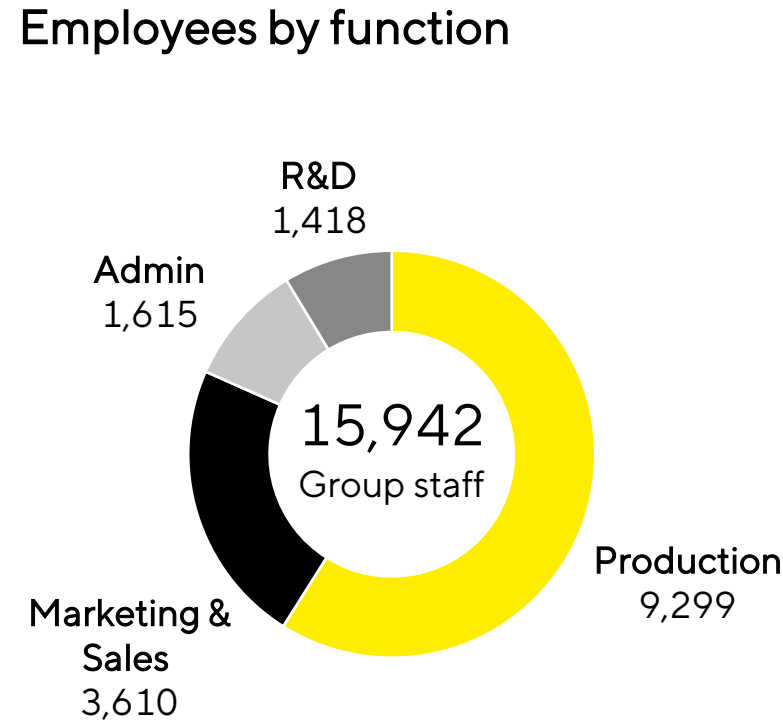
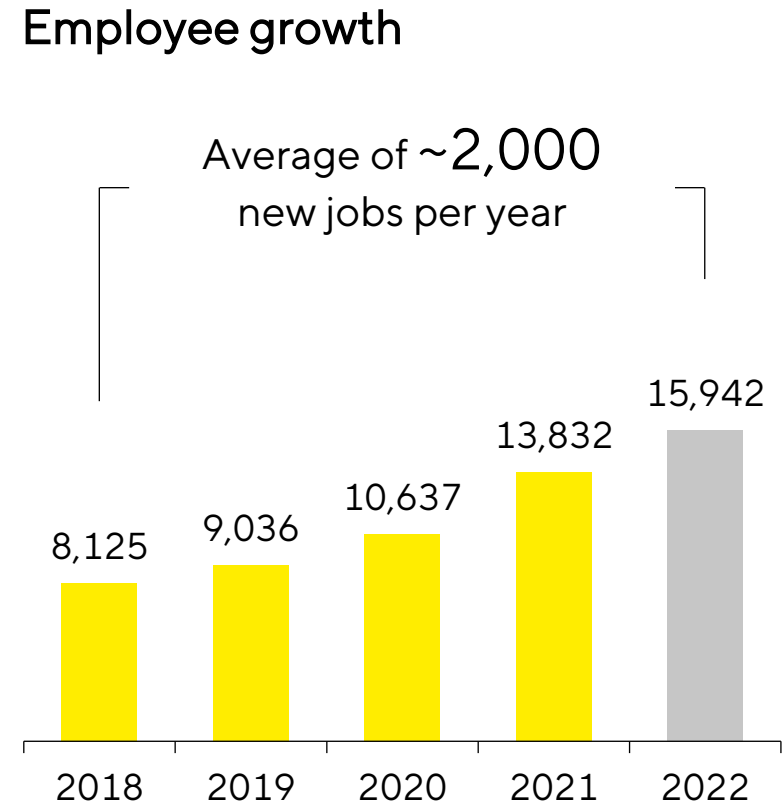
Joachim Kreuzburg, CEO

---



# Welcome 3,897 New People Joined Sartorius in 2022

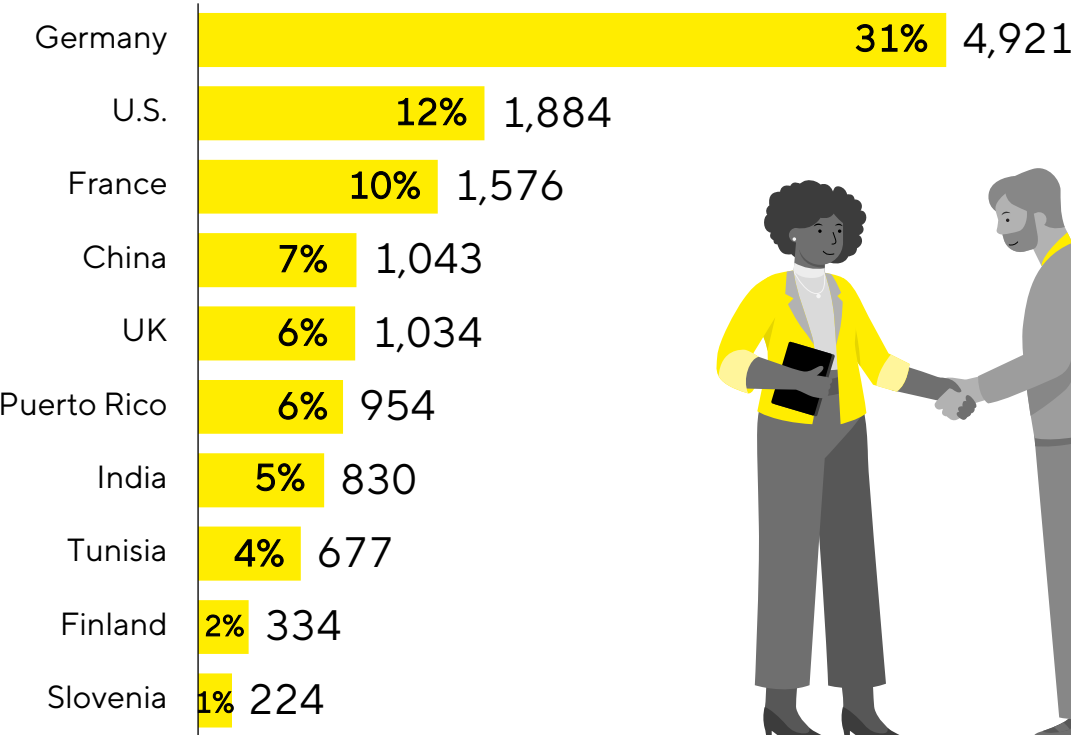
- 3,652 | New hires in 2022
- 245 | New employees from M&A in 2022
- 2,593 | Avg. hires annually 2018-2022
- 60% | Employees with higher education



# Nationality Sartorius Unites People from 119 Nations

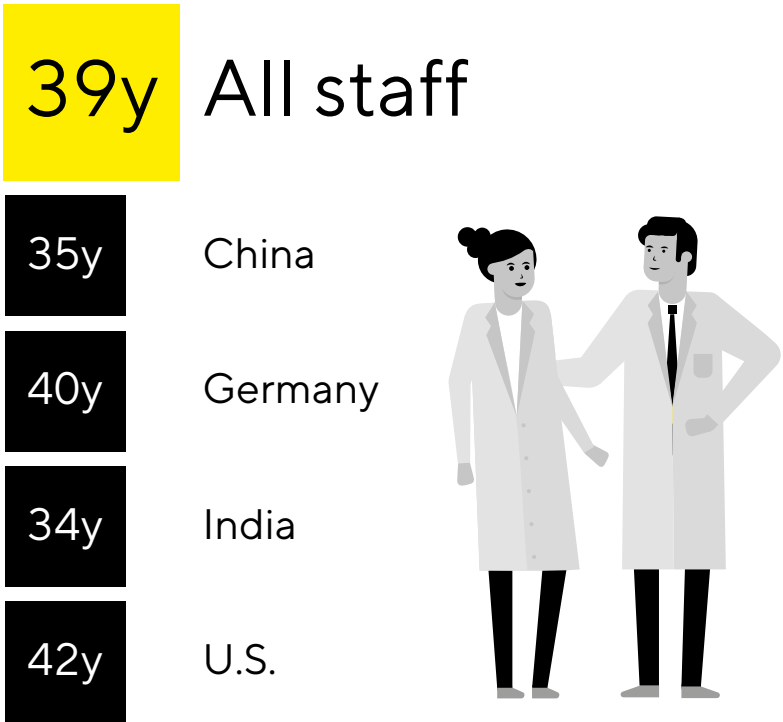


## Top 10 nationalities – Share and number of staff

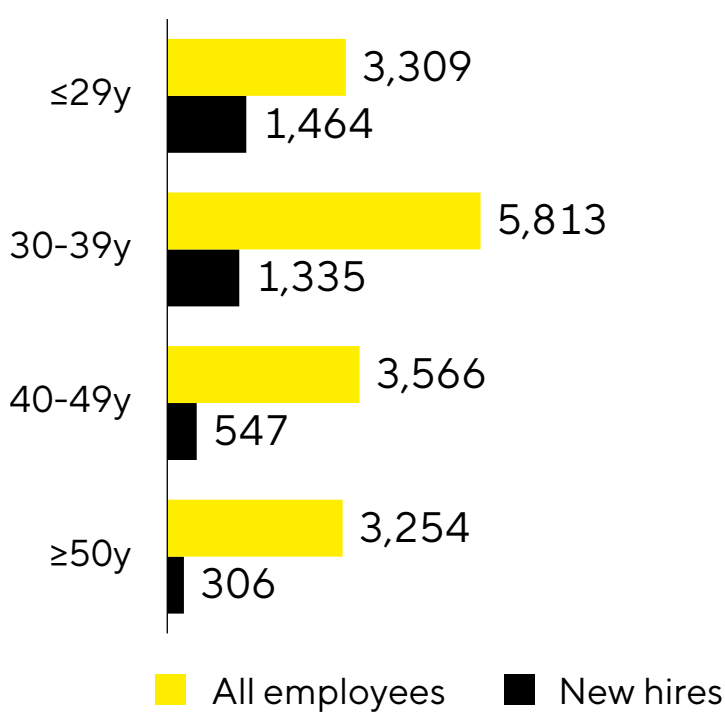


# Age China and India Are Emerging Talent Markets, While Employees 40+ Remain a Valuable Asset

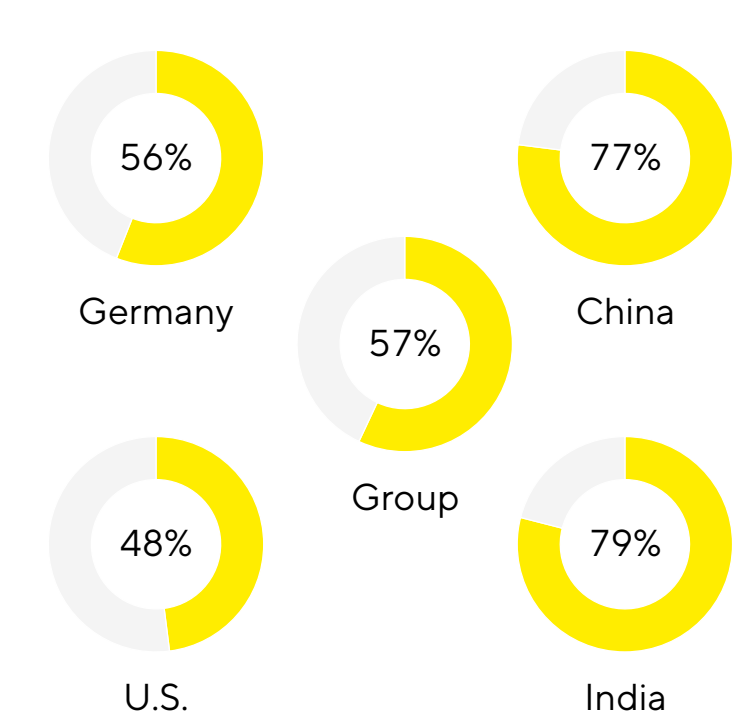
Average age



Number of staff by age



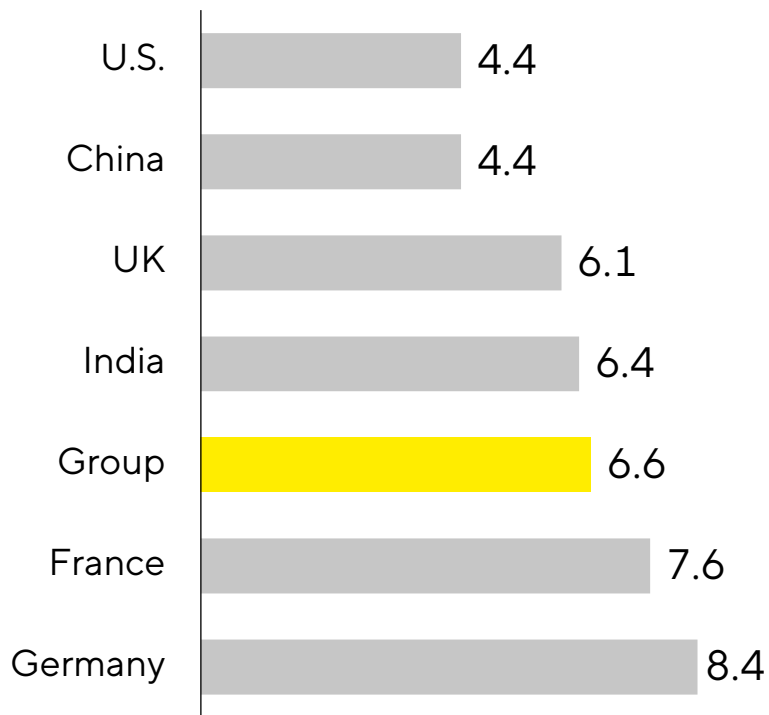
Share of staff younger than 40y



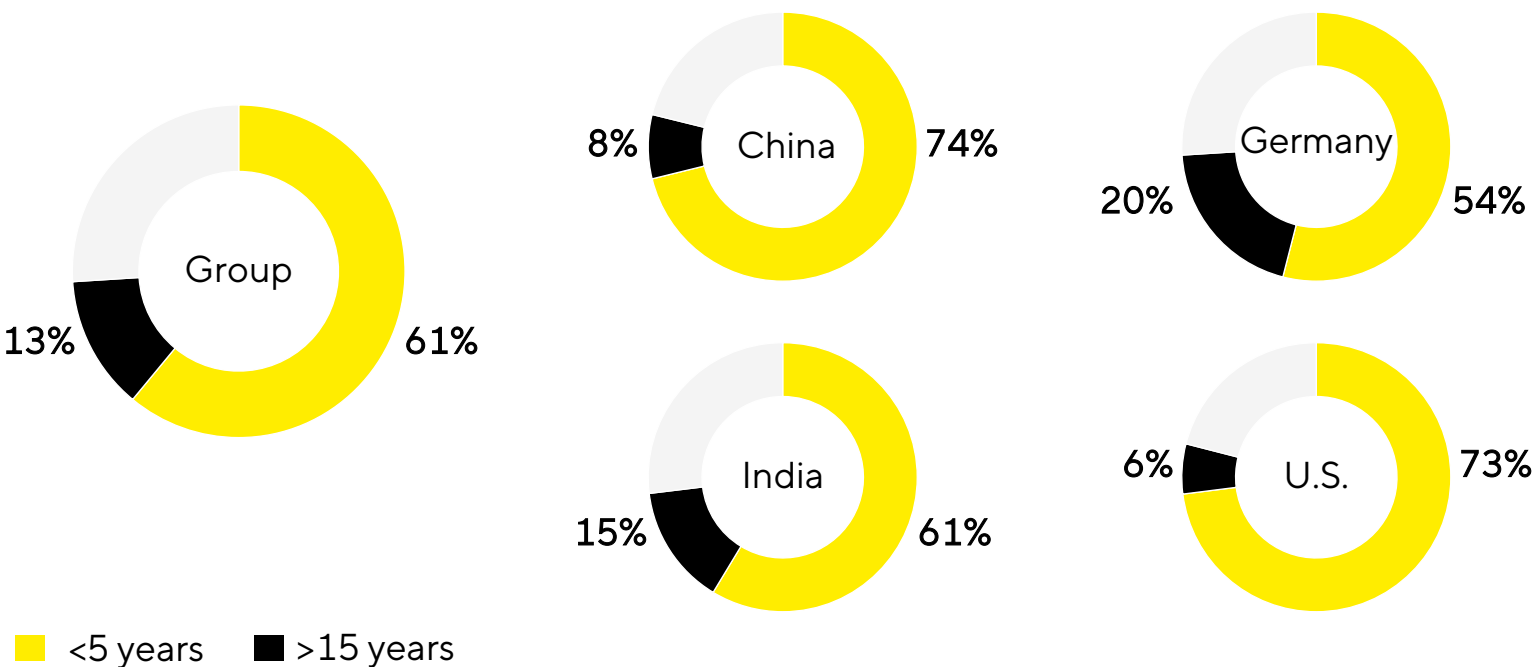


# Tenure Mix of Long-Term Experience and Fresh Ideas

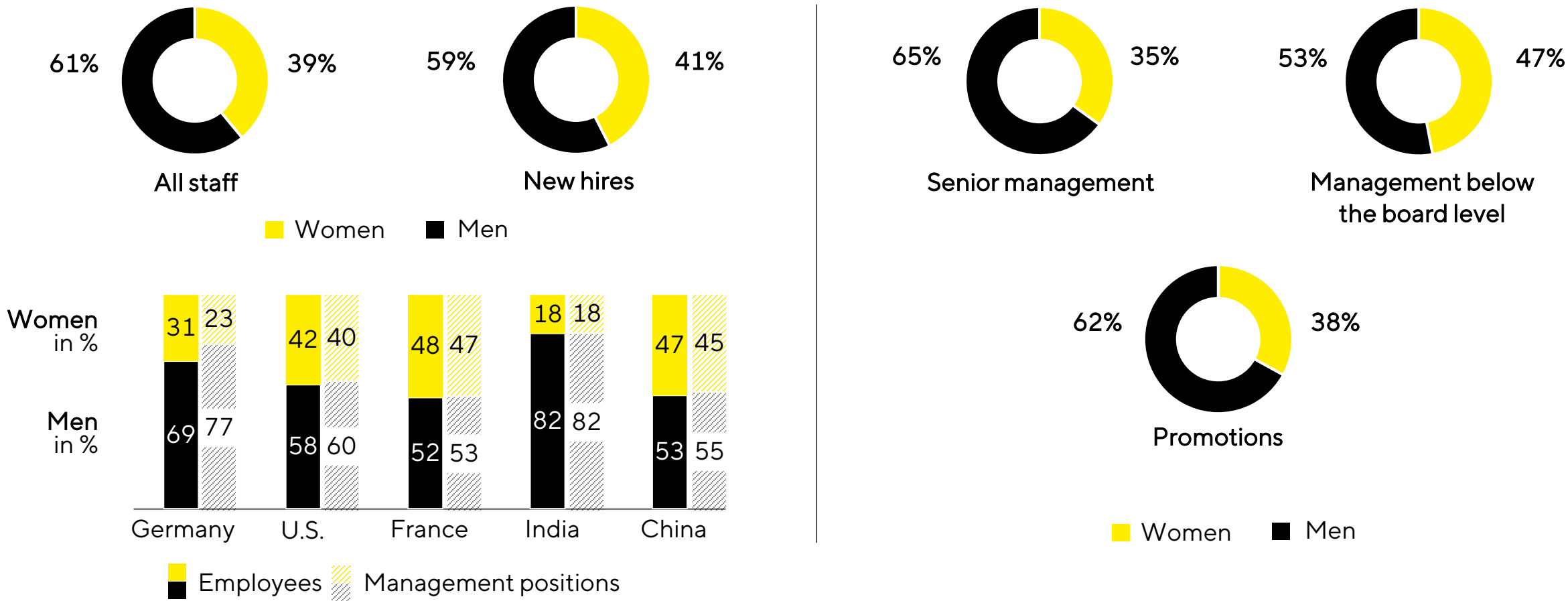
Average tenure in years



Share of staff with tenure higher than 15 years and lower than 5 years



# Gender 47% Women Leaders Below the Board Level



# Fostering exchange and dialogue for women

## **Sartorius Business Women Association**

The network serves to connect women across all regions and aims to support gender parity in leadership positions at Sartorius.

## **Partnership with global non-profit**

Since 2021, Sartorius has been a corporate partner of the Healthcare Businesswomen Association, providing interested employees with access to the organizations network.





# Promoting equality and raising awareness of diversity and inclusion

## North America Diversity, Equity, Inclusion and Belonging (DEIB)

The committee promotes DEIB in the U.S. through community outreach, partnerships, events, and training. So far, 800+ employees have completed Emtram's Unconscious Bias Awareness training.

## LGBT+Friends Alliance

The group strives to enhance the visibility of LGBTQIA+ individuals at Sartorius through virtual events and inclusive "Rainbow Lunches" – creating a safe space for sharing experiences.



# Partnering in the name of diversity and inclusion



Initiative zur Berufs-  
und Studienwahl

# Employee Satisfaction Measured Biannually: 80% Participation Rate



81%  
find that working conditions allow them to fully perform daily work

82%  
experience their team as stable and reliable

84%  
enjoy working for Sartorius

